

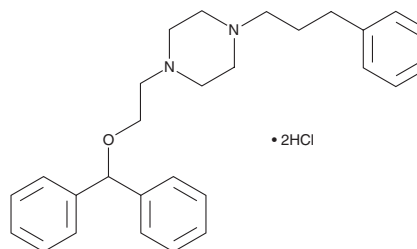
PRODUCT INFORMATION



GBR 12935 (hydrochloride)

Item No. 17700

CAS Registry No.: 67469-81-2
Formal Name: 1-[2-(diphenylmethoxy)ethyl]-4-(3-phenylpropyl)-piperazine, dihydrochloride
MF: C₂₈H₃₄N₂O • 2HCl
FW: 487.5
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

GBR 12935 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the GBR 12935 (hydrochloride) in the solvent of choice. GBR 12935 (hydrochloride) is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of GBR 12935 (hydrochloride) in these solvents is approximately 2 mg/ml.

Description

GBR 12935 is an inhibitor of dopamine uptake that binds to the dopamine transporter ($K_d = 27.2$ nM).^{1,2} It is selective for the dopamine transporter over the serotonin and norepinephrine transporters (K_d s = 940 and 310 nM, respectively).¹ In mice, it increases locomotor activity, but not stereotypic behavior, when administered at a dose of 10 mg/kg but induces gnawing at levels similar to other indirect dopamine agonists with an ED₅₀ value of 17.1 mg/kg.^{2,3}

References

1. Tatsumi, M., Groshan, K., Blakely, R.D., *et al.* Pharmacological profile of antidepressants and related compounds at human monoamine transporters. *Eur. J. Pharmacol.* **340(2-3)**, 249-258 (1997).
2. Tirelli, E. and Witkin, J.M. Differential effects of direct and indirect dopamine agonists on the induction of gnawing in C57Bl/6J mice. *J. Pharmacol. Exp. Ther.* **273(1)**, 7-15 (1995).
3. Tolliver, B.K. and Carney, J.M. Comparison of cocaine and GBR 12935: Effects on locomotor activity and stereotypy in two inbred mouse strains. *Pharmacol. Biochem. Behav.* **48(3)**, 733-739 (1994).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/09/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM